

PRODUCT INFORMATION



FDW028

Item No. 39543

CAS Registry No.: 2768426-49-7

Formal Name: 2-[[[4-ethylphenyl)methyl]amino]-5-[[[(phenylmethyl)amino]methyl]-[1,2,4]triazolo[1,5-a]pyrimidin-7-ol

MF: C₂₂H₂₄N₆O

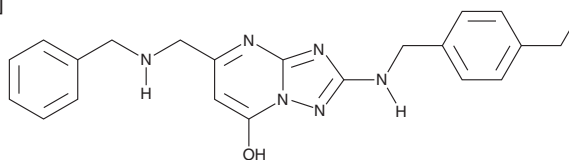
FW: 388.5

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

FDW028 is supplied as a solid. A stock solution may be made by dissolving the FDW028 in the solvent of choice, which should be purged with an inert gas. FDW028 is soluble (≥10 mg/ml) in DMSO.

Description

FDW028 is an inhibitor of fucosyltransferase 8 (FUC8).¹ It binds to FUC8 ($K_d = 5.486 \mu\text{M}$) and inhibits core fucosylation in HCT-8 and SW480 colorectal cancer (CRC) cells. FDW028 (50 μM) induces degradation of B7-H3, also known as CD276, via chaperone-mediated autophagy (CMA) in the same cells. *In vivo*, FDW028 (10 and 20 mg/kg) reduces tumor volume and increases survival in an SW480 mouse xenograft model.

Reference

1. Wang, M., Zhang, Z., Chen, M., *et al.* FDW028, a novel FUT8 inhibitor, impels lysosomal proteolysis of B7-H3 via chaperone-mediated autophagy pathway and exhibits potent efficacy against metastatic colorectal cancer. *Cell Death Dis.* **14**(8), 495 (2023).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/30/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM